

Product Number:	104071
Lot Number:	218805
Expiration:	7-Jul-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Result	pH Target
pH	Pass	7.35 - 7.45

Review/Approval

Sonia Khillan

Name

QC manager

Title

Sonia Khillan / 27 JAN 2023

Signature/Date